## Proteasome 20S α4 subunit monoclonal antibody (MCP79)

This antibody is covered by our Worry-Free Guarantee.

Citations: 5

View Online »

**Ordering Information** 

Order Online »

BML-PW9140-0100

100µl

Manuals, SDS & CofA

**View Online** »

## **Handling & Storage**

**Handling** Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

**Alternative Name** Proteasome subunit α type-7, Proteasome subunit RC6-1,

Proteasome subunit XAPC7

**Application** IHC, WB

**Application Notes** Does not work in IP.

Clone MCP79

**Formulation** Liquid. Ascites in TBS, containing 5 mM sodium azide.

Gene/Protein Identifier PSMA7 (gene name)

**Host** Mouse

**Immunogen** Human placenta derived proteasomes.

lsotype IgG1

**Source** Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Mouse

**Specificity** Recognizes the α4 subunit of the 20S proteasome.

**Technical Info / Product Notes**Various systems for the nomenclature of the proteasome

subunits have been established. This may be a source of confusion as the system on UniProt differs from "standard" nomenclature as described in the literature. The UniProt ID and Gene Name will help to clearly identify the proteins.

014818

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com